

*G-Biosciences/Genotech*: St. Louis, MO, USA; [www.GBiosciences.com](http://www.GBiosciences.com), email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

**Product:** GST Elution Buffer, pH 8.0      **Catalog #:** 786-541      [LU 5.08.09/JK]

<u>Components</u>	<u>Catalog No.</u>	<u>Quantity</u>
GST Elution Buffer, pH 8.0	786-541	100 ml

## **Material Safety Data Sheet**

### **Section 1: Chemical Identification**

**Name:** GST Elution Buffer, pH 8.0

**Catalog #:** a component of product Cat # 786-541 as indicated above

### **Section 2: Composition/Information on Ingredients**

Contains: Glutathione [CAS# 70-18-8] <5%; Tris Base [CAS# 77-86-1] <5%

### **Section 3: Hazards Identification**

Label precautionary statements. Irritating to eyes, respiratory system and skin.

### **Section 4: First-Aid Measures**

**After eye contact:** Immediately flush eyes with copious amounts of water for at least 15 minutes. Call a physician.

**After skin contact:** Immediately wash skin with soap and copious amounts of water.

**After Inhalation:** Remove to fresh air. If not breathing give artificial respiration. If breathing is difficult, call a physician.

**After swallowing:** Wash out mouth with water provided person is conscious. Call a physician.

### **Section 5: Fire Fighting Measures**

**Extinguishing Agents:** Water spray, carbon dioxide, dry chemical powder or appropriate foam.

**Special fire fighting procedures:** None needed.

**Unusual fire and explosions hazards:** None needed.

### **Section 6: Accidental Release Measures**

**Human Safety Precautions:** Wear chemical safety goggles, rubber boots and heavy rubber gloves.

**Cleaning/Collecting Measures:** Mop up, place in a closed container and hold for waste disposal. Ventilate area and wash spill site after material pickup is complete.

### **Section 7: Handling and Storage**

Refer to section 8.

### **Section 8: Exposure Controls/Personal Protection**

**Personal protective equipment:** Chemical safety goggles, compatible chemical-respirator, NIOSH/MSHA-approved respirator, and safety shower and eye bath.

**General Protective and Hygiene Measures:** Mechanical exhaust required.

Avoid contact with eyes, skin and clothing. Wash thoroughly after handling. Keep tightly closed. Store in cool dry place.

**Section 9: Physical and Chemical Properties**

**Form:** Liquid

**Color:** None

**Odor:** None

**Section 10: Stability and Reactivity**

**Thermal Decomposition/Conditions to be Avoided:** No decomposition if used according to specifications.

**Dangerous Reactions:** No dangerous reactions known.

**Dangerous Products of Decomposition:** No known dangerous decomposition products.

**Section 11: Toxicological Information**

**On the skin:** May be harmful if absorbed through the skin. May cause skin irritation.

**On the eye:** May cause eye irritation.

**Inhalation:** May be harmful if inhaled. Material may be irritating to mucous membranes and upper respiratory tract.

**Ingestion:** May be harmful if swallowed.

**Section 12: Ecological Information**

Data not yet available

**Section 13: Disposal Considerations**

**Recommendation:**

Observe all federal, state and local environmental regulations.

**Section 14: Transport Information**

Not dangerous goods. Safe for all transport.

**Section 15: Regulatory information**

Irritating to eyes, respiratory system and skin. In case of contact with eyes, rinse immediately with plenty of water and seek medical advice. Wear suitable protective clothing.

**Section 16: Other Information**

The above information is believed to be correct but does not purport to be all-inclusive and shall be used only as a guide. Geno Technology shall not be held liable for any damage resulting from handling or from contact with the above product. See reverse side of invoice or packing slip for additional terms and conditions of sale.

***G-Biosciences/Genotech St. Louis, MO, USA***

***Web: [www.GBiosciences.com](http://www.GBiosciences.com), email: [technical@GBiosciences.com](mailto:technical@GBiosciences.com)***